

## miProfile™ Human Squamous Cell Carcinoma Exosome miRNA qPCR Array

For focused group profiling of human squamous cell carcinoma exosome related miRNA expression

Cat. No. QM056-A (1 x 96-well plate, Format A)  
 Cat. No. QM056-B (1 x 96-well plate, Format B)  
 Cat. No. QM056-C (1 x 96-well plate, Format C)  
 Cat. No. QM056-D (1 x 96-well plate, Format D)  
 Cat. No. QM056-E (1 x 96-well plate, Format E)

Available as 1 set or 6 sets. Each set contains 37 unique miRNA primers deposited in one 96-well plates

### Introduction

The miProfile™ human squamous cell carcinoma exosome miRNA PCR array is a set of one 96-well plates, covering 37 miRNA primers related to human squamous cell carcinoma exosome. Each 96-well plate contains up to 37 pairs of PCR primers (forward: miRNA-specific primer; reverse: universal primer), which are pre-deposited in each well. Each plate also designated 12 control wells for monitoring the efficiency of every step of the experiment - from reverse transcription to qPCR reaction.

QM056 plate 01: 37 unique miRNA PCR primer pairs

### Shipping and storage condition

Shipped at room temperature  
 Stable for at least 6 months when stored at -20°C

### Array format

GeneCopia provides five qPCR array formats (A, B, C, D, and E) suitable for use with the following real-time cyclers.

**Important note:** Upon receiving, please check to make sure that the correct array format was ordered to ensure the compatibility with your qPCR instrument.

| Plate format          | Instrument provider   | qPCR instrument model  |
|-----------------------|-----------------------|--|
| <b>A</b><br>(96-well) | Applied Biosystems    | 5700, 7000, 7300, 7500, 7700, 7900HT (Standard 96-well block), ViiA™7 (Standard 96-well block) |
| <b>B</b><br>(96-well) | Applied Biosystems    | 7500 (Fast block), 7900HT (Fast block), StepOnePlus™, ViiA™7 (Fast block)                      |
| <b>C</b><br>(96-well) | Bio-Rad Laboratories  | iCycler iQ®, MyiQ™, iQ™5   |
| <b>D</b><br>(96-well) | Bio-Rad Laboratories  | CFX96™, DNA Engine Opticon™, DNA Engine Opticon 2™, Chromo4™                                   |
| <b>E</b><br>(96-well) | Roche Applied Science | LightCycler® 480 (96-well block)   |

## Array layout

|   | 1  | 2  | 3   | 4   | 5   | 6   | 7   | 8   | 9  | 10 | 11  | 12  |
|---|----|----|-----|-----|-----|-----|-----|-----|----|----|-----|-----|
| A | 1  | 2  | 3   | 4   | 5   | 6   | 7   | 8   | 9  | 10 | 11  | 12  |
| B | 13 | 14 | 15  | 16  | 17  | 18  | 19  | 20  | 21 | 22 | 23  | 24  |
| C | 25 | 26 | 27  | 28  | 29  | 30  | 31  | 32  | 33 | 34 | 35  | 36  |
| D | 37 | 38 | 39  | 40  | 41  | 42  | 43  | 44  | 45 | 46 | 47  | 48  |
| E | 49 | 50 | 51  | 52  | 53  | 54  | 55  | 56  | 57 | 58 | 59  | 60  |
| F | 61 | 62 | 63  | 64  | 65  | 66  | 67  | 68  | 69 | 70 | 71  | 72  |
| G | 73 | 74 | 75  | 76  | 77  | 78  | 79  | 80  | 81 | 82 | 83  | 84  |
| H | NC | NC | HK1 | HK2 | HK3 | HK4 | HK5 | HK6 | RT | RT | PCR | PCR |

**Figure 1. Illustration of miProfile miRNA qPCR array layout (96-well plate)**

**miRNA primer pairs:** Wells 1-84 are designated wells for pre-deposited miRNA primer pairs.

**NC:** Negative controls, which only have the pre-deposited reverse universal primers

**HK1-6:** Six pre-deposited housekeeping snRNAs primer pairs, which can be used as endogenous positive controls as well as for array normalization.

**RT:** Two replicates of spike-in reverse transcription controls, which can be used to monitor the efficiency of the RT reaction. These pre-deposited primer pairs specifically amplify the cDNA template reversed transcribed from the spike-in exogenous RNA in the sample.

**PCR:** Two replicates of positive PCR controls, which are used to verify the PCR efficiency by amplifying the pre-deposited DNA template with its specific pre-deposited primer pairs.

## Materials required but not provided

All-in-One™ miRNA First-Strand cDNA Synthesis Kit  
 All-in-One™ qPCR Mix  
 Total RNA extraction kit (RNAzol® RT RNA extraction reagent is recommended)  
 DNase/RNase free tips, PCR reaction tubes, 1.5 ml microcentrifuge tubes  
 5 ml and 10 ml graduated pipettes, beakers, flasks, and cylinders  
 10 µl to 1,000 µl adjustable single channel micropipettes with disposable tips  
 5 µl to 20 µl adjustable multichannel micropipette, disposable tips, and reservoir  
 qPCR instrument, compatible with miRNA qPCR arrays ordered

## miRNA primer list

The primer list can be downloaded from <http://www.genecopoeia.com/product/qpcr-arrays/mirna/cancer.php>

## Limited Use License

Following terms and conditions apply to use of miProfile Human Brain Cancer miRNA qPCR Arrays (the Product). If the terms and conditions are not acceptable, the Product in its entirety must be returned to GeneCopoeia within 5 calendar days. A limited End-User license is granted to the purchaser of the Product. The Product shall be used by the purchaser for internal research purposes only. The Product is expressly not designed, intended, or warranted for use in humans or for therapeutic or diagnostic use. The Product must not be resold, repackaged or modified for resale, or used to manufacture commercial products or deliver information obtained in service without prior written consent from GeneCopoeia. This Product should be used in accordance with the NIH guidelines developed for recombinant DNA and genetic research. Use of any part of the Product constitutes acceptance of the above terms.

### Limited Warranty

GeneCopoeia warrants that the Product meets the specifications described in the accompanying Product Datasheet. If it is proven to the satisfaction of GeneCopoeia that the Product fails to meet these specifications, GeneCopoeia will replace the Product. In the event a replacement cannot be provided, GeneCopoeia will provide the purchaser with a refund. This limited warranty shall not extend to anyone other than the original purchaser of the Product. Notice of nonconforming products must be made to GeneCopoeia within 30 days of receipt of the Product. GeneCopoeia's liability is expressly limited to replacement of Product or a refund limited to the actual purchase price. GeneCopoeia's liability does not extend to any damages arising from use or improper use of the Product, or losses associated with the use of additional materials or reagents. This limited warranty is the sole and exclusive warranty. GeneCopoeia does not provide any other warranties of any kind, expressed or implied, including the merchantability or fitness of the Product for a particular purpose.

GeneCopoeia is committed to providing our customers with high-quality products. If you should have any questions or concerns about any GeneCopoeia products, please contact us at 301-762-0888.

© 2019 GeneCopoeia, Inc.

GeneCopoeia, Inc.  
9620 Medical Center Drive, Suite 101  
Rockville, MD 20850  
+1 (301) 762-0888  
+1 (866) 360-9531  
[inquiry@genecopoeia.com](mailto:inquiry@genecopoeia.com)